

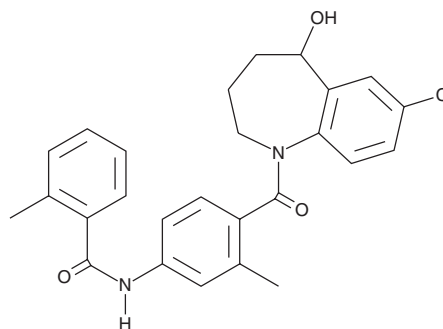
PRODUCT INFORMATION



Tolvaptan

Item No. 19691

CAS Registry No.: 150683-30-0
Formal Name: N-[4-[(7-chloro-2,3,4,5-tetrahydro-5-hydroxy-1H-1-benzazepin-1-yl)carbonyl]-3-methylphenyl]-2-methyl-benzamide
Synonym: OPC 41061
MF: C₂₆H₂₅ClN₂O₃
FW: 448.9
Purity: ≥98%
UV/Vis.: λ_{max}: 269 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Tolvaptan is supplied as a crystalline solid. A stock solution may be made by dissolving the tolvaptan in the solvent of choice, which should be purged with an inert gas. Tolvaptan is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of tolvaptan in ethanol is approximately 12 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Tolvaptan is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tolvaptan should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Tolvaptan has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Tolvaptan is a nonpeptide vasopressin V₂ receptor antagonist (IC₅₀ = 3 nM for rat receptor) and a diuretic agent.¹ It is selective for V₂ over V₁ receptors (IC₅₀ = 0.58 μM). Tolvaptan increases urine volume by 3-fold in rats when administered at a dose of 0.54 mg/kg. It also reduces left ventricular end-systolic volumes and improves left ventricular ejection fraction in a rat model of myocardial infarction.^{43556} Formulations containing tolvaptan have been used to treat hyponatremia.

Reference

1. Kondo, K., Ogawa, H., Yamashita, H., *et al.* 7-Chloro-5-hydroxy-1-[2-methyl-4-(2-methylbenzoyl-amino) benzoyl]-2,3,4,5-tetrahydro-1H-1-benzazepine (OPC-41061): A potent, orally active nonpeptide arginine vasopressin V₂ receptor antagonist. *Bioorg. Med. Chem.* **7(8)**, 1743-1754 (1999).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/14/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM